

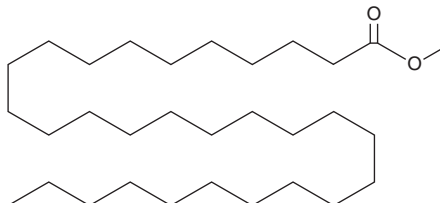
# PRODUCT INFORMATION



## Triacontanoic Acid methyl ester

Item No. 26727

**CAS Registry No.:** 629-83-4  
**Formal Name:** triacontanoic acid, methyl ester  
**Synonyms:** C30:0 methyl ester, Methyl Melissate, Methyl Triacontanoate, NSC 20743, SFE 31:0  
**MF:**  $C_{31}H_{62}O_2$   
**FW:** 466.8  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Triacontanoic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the triacontanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Triacontanoic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of triacontanoic acid methyl ester in ethanol and DMF is approximately 25 mg/ml and approximately 10 mg/ml in DMSO.

Triacontanoic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triacontanoic acid methyl ester should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Triacontanoic acid methyl ester has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Triacontanoic acid methyl ester is a fatty acid methyl ester that has been found in sediment samples from the Harney River in Florida and Lake Kivu in the East African rift valley.<sup>1,2</sup>

### References

1. Jaffé, R., Rushdi, A.I., Medeiros, P.M., *et al.* Natural product biomarkers as indicators of sources and transport of sedimentary organic matter in a subtropical river. *Chemosphere* **64**(11), 1870-1884 (2006).
2. Al-Mutlaq, K., Standley, L.J., and Simoneit, B.R. Composition and sources of extractable organic matter from a sediment core in Lake Kivu, East African rift valley. *Appl. Geochem.* **23**(5), 1023-1040 (2008).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/13/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM